



## **MyBath™ 4 Mini Water Bath - Model B2000-4 Instruction Manual**

Benchmark's MyBath 4 is a micro-processor controlled - thermostatic water bath and is widely used for petroleum, biochemical and scientific research. The instrument is designed for accurate temperature control between ambient +5°C up to 99.9°C.

### **I. Product Specifications:**

**Capacity:** 4L

**Temperature Range:** Ambient +5°C to 99.9°C.

**Temperature Accuracy:** +/-0.2°C (37°C)    **Temperature Uniformity:** +/-0.2°C (37°C)

**Dimensions (Internal):** wxdxh 13.5 x 24 x 14 cm (5.3 x 9.4 x 5.5")

**Dimensions (External):** wxdxh 19.5 x 36 x 24 cm (7.5 x 14 x 9.5")

**Weight:** 6.5 kg (15 lbs)    **Electrical:** 100-120VAC or 220-240VAC, 50/60Hz, 8A/5A

**Operating Environment:** 10°C to 30°C    70%RH    **Warranty:** 1 year

### **II. Product Set-Up:**

Place the water bath on a clean, level and stable surface. Make sure that the power switch is in the OFF position. Plug the power cord into the back of the unit, then plug the electrical cord into a properly grounded outlet of the correct voltage (shown on the serial number label). Prior to powering up the instrument, fill the tank with water (at least 50%).

**WARNING:** Before heating, ALWAYS ensure that the water tank is at least 50% full.

### **III. Product Operation:**

Once the tank is filled with liquid, power on the instrument (power switch is located on the back of the instrument).

To set the desired temperature, press the "Temp." button and the current set temperature will flash on the display. Now use the up and down arrows to set the desired temperature.

When the instrument is in the heating state, the hot indication light will illuminate.

**WARNING:** NEVER touch the inner chamber when the hot indication light is illuminated.

### **IV. Recalibration:**

If adjustment to the temperature calibration is required, this can be accomplished by following the steps below:

- Set the desired temperature to perform the calibration
- After this desired temperature has been reached, insert a thermometer and record the temperature
- Press the "Cal." Key and the displayed temperature flashes. Now use the up and down arrows to set the recorded temperature on the display. After 8 seconds the calibration is stored.

### **V. Care and Maintenance:**

No routine maintenance is required other than to keep the unit clean. Cleaning can be done with a damp cloth. Avoid the use of solvents as they may attack the product housing.